



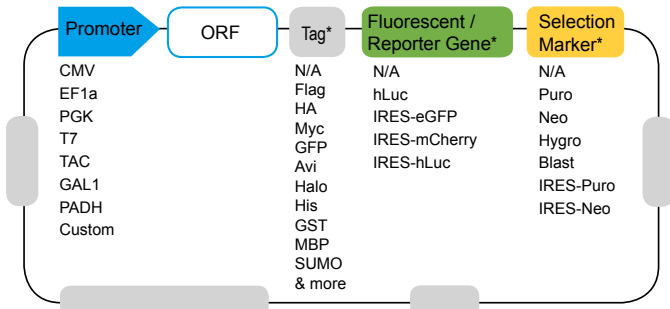
# OmicsLink™ Expression-Ready ORF Clones

## Nextday-ORF™ Collection

140,000 clones Starting from \$69

Vector	Promoter	Vector Type	Marker	Tag
M02	CMV	Mammalian	Neo	None
M13	CMV	Mammalian	Neo	C-Flag
M98	CMV	Mammalian	Neo	C-eGFP
Lv105	CMV	Lentiviral	Puro	None
Lv242	CMV	Lentiviral	Puro	C-Avi-Flag

## OmicsLink™ ORF Clone Collections



\* Visit our website for a complete list of tags, fluorescent/reporter genes and selection markers.

## Advantages

- Starting from \$69 for selected clones
- Next-day shipping of 140,000 Nextday-ORF™ clones
- All ORFs from human and mouse **sequence validated**
- 200+ expression vectors for mammalian, lentiviral and more



## Clone Sets

Human Gene or Custom Design

Set type	# of Genes	# of Clones
CDs	341	430
Cytokines and cytokine receptors	347	420
Drug target genes	5022	6144
GPCRs	526	583
Innate immune proteins	565	702
Ion channels	297	366
Kinases	707	940
Nuclear hormone receptors	56	82
Phosphatases	213	261
Protease inhibitors	111	134
Secretory proteins	1497	1788

For more clone sets, visit our website [www.genecopoeia.com](http://www.genecopoeia.com).

## Related Products

- Custom gene synthesis
- Custom Cloning
- Lentiviral & AAV particles
- CRISPR-Cas9 Services
- EZShuttle™ Gateway® PLUS ORF Clones

## Contact us:

### GeneCopoeia, Inc.

9620 Medical Center Dr. Suite 101,  
Rockville, MD 20850 USA



Email	<a href="mailto:inquiry@genecopoeia.com">inquiry@genecopoeia.com</a>
Tel	+1 (301) 762-0888
Toll free	+1 (866) 360-9531
Fax	+1 (301) 762-3888
Website	<a href="http://www.genecopoeia.com">www.genecopoeia.com</a>

# OmicsLink™ Expression-Ready ORF Clones

## Selected Mammalian Expression Vectors

Vector	promoter	Tag/Reporter	Selection Marker
M02	CMV	-	Neo
M13	CMV	C-Flag	Neo
M98	CMV	C-eGFP (monomeric)	Neo
M61	CMV	IRES2-eGFP	Neo
M56	CMV	C-mCherry	Neo
M29	CMV	N-eGFP	Neo
M03	CMV	C-eGFP	Neo
M06	CMV	N-3×HA	Neo
M11	CMV	N-Flag	Neo
M09	CMV	C-Myc	Neo
M12	CMV	N-3×Flag	Neo
M55	CMV	N-mCherry	Neo
M14	CMV	C-3×Flag	Neo
M07	CMV	C-3×HA	Neo
M77	CMV	C-His	Neo
M01	CMV	N-His	Neo
M43	CMV	N-Myc	Neo
M46	CMV	C-Flag-IRES-eGFP	Neo
M51	CMV	C-His-IRES-eGFP	Neo
M16	CMV	C-eYFP	Neo
M10	CMV	C-Myc-His	Neo
M45	CMV	C-3×HA-IRES-eGFP	Neo
M72	CMV	C-Myc-IRES2-eGFP	Neo
M15	CMV	N-eYFP	Neo
M17	CMV	C-Avi	Neo
M48	CMV	N-Avi-IRES-Biotin ligase	Neo
M50	CMV	C-Halo/Tev protease	Neo
M08	CMV	C-3×HA-His	Neo
M49	CMV	N-Halo/Tev protease	Neo
M05	CMV	N-Avi	Neo
M04	CMV	N-GST/EK protease	Neo
M33	CMV	C-eCFP	Neo
M32	CMV	N-eCFP	Neo

## Selected Lentiviral Expression Vectors

Vector	promoter	Tag/Reporter	Selection Marker
Lv105	CMV	-	Puro
Lv152	CMV	-	Hygro
Lv242	CMV	C-Avi-Flag	Puro
Lv205	CMV	IRES2-eGFP	IRES-Puro
Lv120	CMV	C-3×HA	Puro
Lv102	CMV	N-Flag	Puro
Lv203	CMV	C-Flag-SV40-eGFP	IRES-Puro
Lv121	CMV	C-Flag	Puro
Lv201	CMV	SV40-eGFP	IRES-Puro
Lv207	CMV	SV40-eGFP	IRES-Hygro
Lv130	CMV	C-mCherry	Puro
Lv213	CMV	IRES2-mCherry	IRES-Puro
Lv103	CMV	N-eGFP	Puro
Lv216	CMV	C-Flag-SV40-mCherry	Puro
Lv125	CMV	C-Myc	Puro
Lv111	CMV	N-mCherry	Puro
Lv217	CMV	SV40-hLUC	IRES-Puro
Lv129	CMV	C-Halo	Puro
Lv211	CMV	C-Halo	IRES-Hygro
Lv206	CMV	SV40-mCherry	IRES-Puro
Lv109	CMV	N-eCFP	Puro
Lv128	CMV	C-His	Puro
Lv110	CMV	N-Halo	Puro
Lv123	CMV	C-eYFP	Puro
Lv126	CMV	C-Myc-His	Puro
Lv202	CMV	C-Myc-SV40-eGFP	IRES-Puro
Lv204	CMV	C-Myc-SV40-mCherry	IRES-Puro
Lv104	CMV	N-eYFP	Puro
Lv108	CMV	N-Avi	Puro
Lv127	CMV	C-eCFP	Puro

Not all available tags, reporter genes and selection markers are listed here. Please visit our website [www.genecopoeia.com](http://www.genecopoeia.com) for a complete list .